



## JBScreen Basic 1 - 4

Cat. No.	Amount
CS-125	4 Kits

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at 8-10 °C

**Shelf Life:** 12 months